



---

## **PRODUCT INFORMATION**

Product  VECTASHIELD® Mounting Medium with PI

Catalog No.  H-1300

Amount  10 ml

Concentration of PI  1.5 µg/ml

Storage Conditions  Refrigerate in the dark

This lot has been tested for photobleaching retardance with qualitative and quantitative methods using fluorescent microscopy.

VECTASHIELD® Mounting Medium is a unique, stable formula for preserving fluorescence. VECTASHIELD® Mounting Medium prevents rapid photobleaching of fluorescein, AMCA, Alexa Fluor® 488 and other fluorochromes.\* Sections mounted in VECTASHIELD® Mounting Medium retain fluorescence during prolonged storage.

VECTASHIELD® Mounting Medium is supplied in a bottle fitted with a cone-seal cap. A cap fitted with a drop dispenser pipet is also included. The cone-seal cap can be replaced with the drop dispenser cap for routine use and during storage. Although the pipet bulb is made of inert materials, it is suggested that once the drop dispenser cap has been secured on the bottle, the bottle should be stored in an upright position.

To mount tissues or cells on a slide, dispense one drop of VECTASHIELD® Mounting Medium onto the section. Then coverslip and allow VECTASHIELD® Mounting Medium to disperse over the entire section. Applying an excess of mounting medium is not recommended. The special pipet supplied with the product is designed to deliver small drop volumes of approximately 25 µl.

\* VECTASHIELD® is not recommended for use with Cy2™, which may dissipate in the section over time.



VECTASHIELD® Mounting Medium does not solidify, but remains a liquid on the slide. After mounting, coverslipped slides will not readily dry out and can be reviewed for weeks afterward without sealing. For prolonged storage coverslips can be permanently sealed around the perimeter with nail polish or a plastic sealant. Mounted slides should be stored at 4 °C protected from light.

This special formulation of VECTASHIELD® Mounting Medium contains propidium iodide (PI) and is intended to be used to counterstain DNA. This product is designed to be used for *in situ* hybridization techniques or for other procedures requiring fluorescent labeling of DNA. PI excites at about 535 nm and emits at about 615 nm when bound to DNA, producing a red fluorescence. PI will also stain RNA. Although PI is a mutagen and appropriate precautions should be taken when using it, the concentration of PI is so low in this product that it is not appropriate to list it in hazard data information.

*VECTASHIELD* is a registered trademark of Vector Laboratories, Inc.  
*Alexa Fluor* is a registered trademark of Molecular Probes